



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP3-00495

NBP3-00495 Protocol

Flow Cytometry Protocol for Absolute Rainbow Cell Count Particle Set (NBP3-00495)

Shake bottle vigorously or vortex briefly before use.

1. Add the monoclonal antibody of your choice to 100 μ L of test sample.
2. Incubate, lyse, wash, and then resuspend in 1 to 2 mL PBS. If staining and lysing are not needed, just mix a known volume of test sample in 1 to 2 mL PBS
3. Add exactly 50 μ L of NBP3-00495 particles to the suspension. Precision when pipetting is absolutely critical.
4. Run the sample in the flow cytometer and obtain the amount of events for the NBP3-00495 particles and your test sample.
5. Calculate the number of cells accordingly:
 $(A/B) \times (C/D) = \text{Number of cells per } \mu\text{L}$
where
A = number of events for the test sample
B = number of events for the NBP3-00495 particles
C = number of NBP3-00495 particles per 50 μ L
D = volume of test sample initially used